

# TaqMan® 3C Chromosome Conformation Kits: 3C Library Reagents

Pub Part Number 4466153 Revision Date 15 June 2012 (Rev. D)

Cat. no.	Box	Contents			Storage conditions
4466151 (EcoRI) 4466152 (HindIII) 4477132 (Library Reagents only)	3C Library Reagents (Ambient)	3C Quenching Solution	3C RE Stop Solution	3C DNA Wash Solution	Room temperature (RT, 18°C to 25°C)
		3C Neutralization Solution	3C Elution Solution 1 (ES1)	3C Elution Solution 2 (ES2)	
		3C 1X PBS	3C DNA Binding Solution		4°C
	3C Library Reagents (-20°C)	3C Lysis Solution	3C Magnetic Beads		
		3C Protease Inhibitor	3C DNA Ligase	3C RNase A	-15°C to -25°C
		3C 5X Ligation Buffer	3C Proteinase K		

**Note:** For safety and biohazard guidelines, refer to the “Safety” section in the *TaqMan® 3C Chromosome Conformation Kits User Guide – Early Access* (Pub no. 4468142). For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Related documentation

The following related documents are available at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support):

Document	Part number	Description
<i>TaqMan® 3C Chromosome Conformation Kits User Guide – Early Access</i>	4468142	Provides detailed procedures for using the 3C Library Reagents and 3C Human Control Assays, guidelines for designing assays, and troubleshooting information.
<i>TaqMan® 3C Chromosome Conformation Kits Quick Reference – Early Access</i>	4471872	Provides brief procedures for advanced users who are familiar with the procedures in the <i>TaqMan® 3C Chromosome Conformation Kits User Guide – Early Access</i> .

## Early Access Disclaimer

Customer acknowledges that this product is not a commercially released product and is still under development. THIS PRODUCT AND ASSOCIATED MATERIALS ARE SUPPLIED "AS IS" WITHOUT ANY WARRANTY, EXPRESS OR IMPLIED, AND ALL WARRANTIES ARE EXPRESSLY DISCLAIMED, INCLUDING WITHOUT LIMITATION THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, WHETHER ARISING FROM A STATUTE OR OTHERWISE IN LAW OR FROM A COURSE OF DEALING OR USAGE OF TRADE. Nothing herein shall be deemed to be or imply a duty for Life Technologies and/or its affiliate(s) to further develop this product or to release a commercial version thereof.



**For Research Use Only. Not intended for human or animal therapeutic or diagnostic use.**

NOTICE TO PURCHASER: PLEASE REFER TO THE TAQMAN® 3C CHROMOSOME CONFORMATION KITS USER GUIDE — EARLY ACCESS FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.

**DISCLAIMERS**

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license.

**For support visit** [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support)

[www.lifetechnologies.com](http://www.lifetechnologies.com)

